



# **Alcohols and acetates**

## Analysis of volatiles in distilled spirits

### Application Note

Food Testing & Agriculture

#### **Authors**

Agilent Technologies, Inc.

#### **Introduction**

Gas chromatography with an Agilent CP-Wax 57 CB for Glycols and Alcohols column separates ten volatiles in distilled spirits in eight minutes.



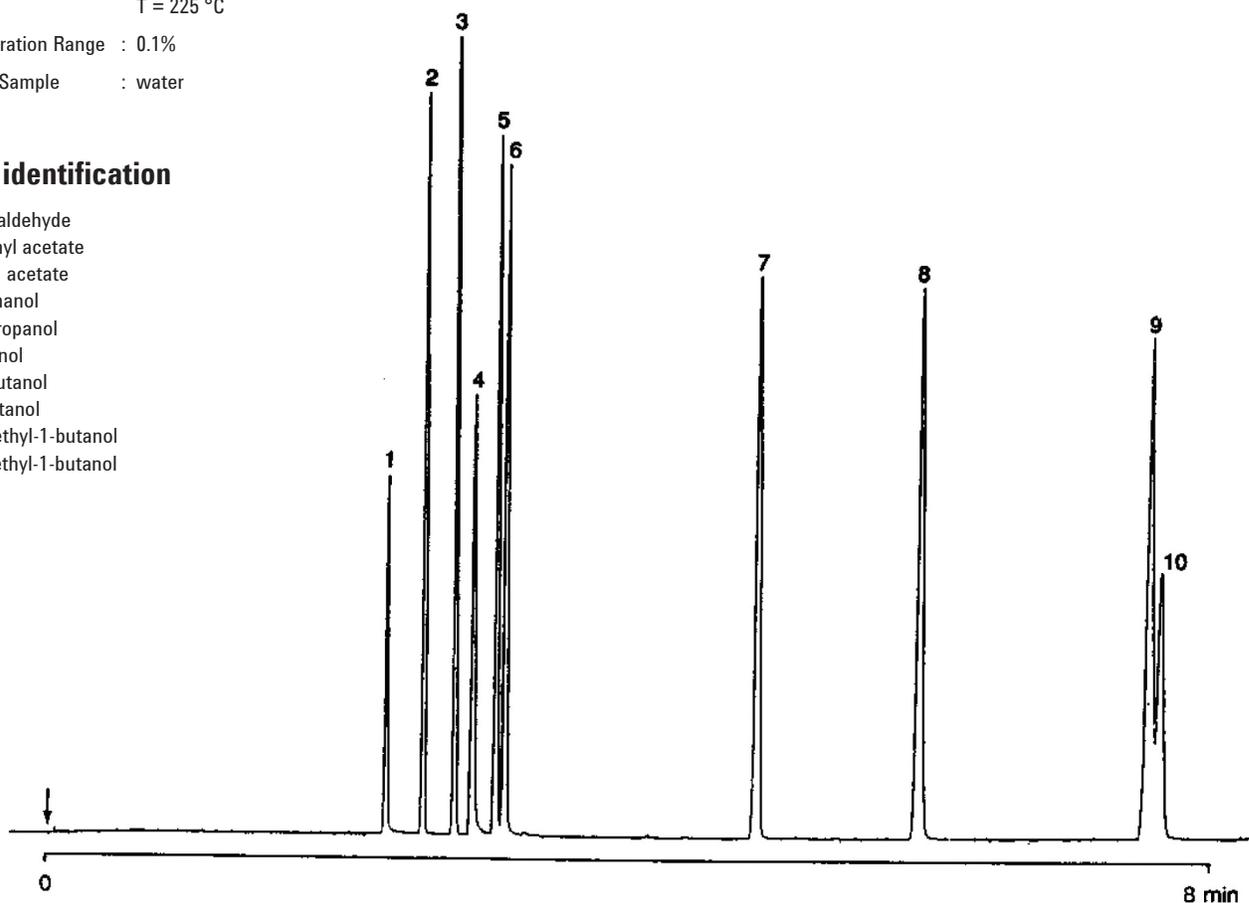
**Agilent Technologies**

## Conditions

Technique : GC-capillary  
Column : Agilent CP-Wax 57 CB for glycols and alcohols, 0.25 mm x 25 m fused silica WCOT CP-Wax 57 CB (df = 0.2  $\mu$ m) (Part no. CP7615)  
Temperature : 60 °C  $\rightarrow$  80 °C, 2 °C/min; 80 °C  $\rightarrow$  150 °C, 10 °C/min  
Carrier Gas : N<sub>2</sub>, 50 kPa (0.5 bar, 7.3 psi), 19 cm/s  
Injector : Splitter, 60 mL/min  
T = 200 °C  
Detector : FID, 4 x 10<sup>-12</sup> Afs  
T = 225 °C  
Concentration Range : 0.1%  
Solvent Sample : water

## Peak identification

1. acetaldehyde
2. methyl acetate
3. ethyl acetate
4. methanol
5. isopropanol
6. ethanol
7. isobutanol
8. n-butanol
9. 2-methyl-1-butanol
10. 3-methyl-1-butanol



[www.agilent.com/chem](http://www.agilent.com/chem)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A00692



**Agilent Technologies**